

Product Number:	103718
Lot Number:	503720
Expiration:	29-Nov-2022



Component	Item Number	Lot Number
SBG Concentrate	103650	227217

SBG Concentrate Concentration (nM)	160
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 13 OCT 2022
 Name Title Signature/Date